

PRODUCT INFORMATION



PSSG Blocking Reagent

Item No. 10010727

Overview

Contents: This vial contains a crystalline solid.
Synonym: Protein S-glutathionylation Blocking Reagent
Stability: ≥3 years at -20°C

Laboratory Procedures

For use in Cayman's S-Glutathionylated Protein Detection Kit (Item No. 10010721). Three vials are provided in the kit to allow for up to three separate experiments. Prior to use, reconstitute the contents of a vial by the addition of 100 µl of DMF to the lyophilized solid, followed by mixing and dilution to 10 ml final volume with Cayman's 1X PSSG Assay Buffer (Item No. 10010726). Do not store any unused reconstituted reagent for more than one day. *Dispose of unused reagent in aqueous hazardous waste.*

Precaution

This product may not perform as described if used with other assay reagents or protocols.

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/23/2023

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM